

## SUV39H2-[CHR] (His)

CATALOG NO.: RD-11-466

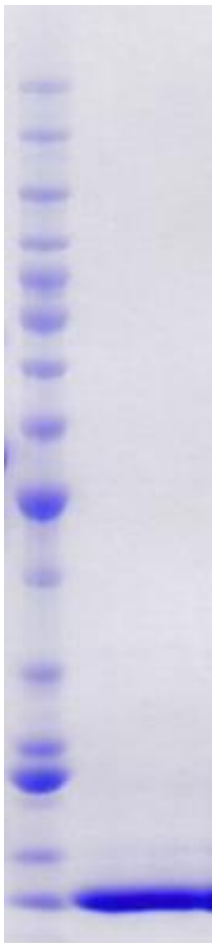
LOT NO.:

**DESCRIPTION:** Human recombinant SUV39H2 chromodomain region (residues 46-108; Genbank Accession # NM\_001193424.1; MW = 10.3 kDa) expressed as an N-terminal His-tag fusion protein in *E. coli*.

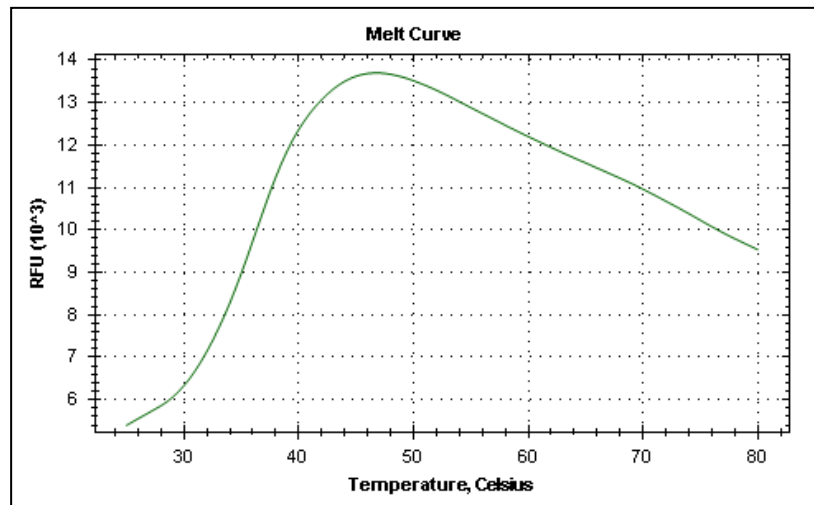
**PURITY:** >90% by SDS-PAGE

**SUPPLIED AS:** \_\_\_ µg/µL in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 10% glycerol, 1mM TCEP as determined by OD<sub>280</sub>.

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry ice/ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 µg of RBC SUV39H2-[CHR] (His). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**Differential Scanning Fluorimetry of RBC SUV39H2-[CHR] (His).** Thermal denaturation of MPP\*-[CHR] (His) is detected (CFX384™ Touch thermal cycler, 'FRET' channel; Bio-Rad) by increased binding and fluorescence of the dye SYPRO® Orange (Life Technologies). The apo form of SUV39H2-[CHR] (His) displays a T<sub>m</sub> of 36.5°C.

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.